



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

RNaseReveal™ Activity Assay Kit

Fluorescent assay for sensitive detection of RNase activity in liquid samples. Ideal for detecting RNase contamination in quality control workflows.



Product attributes

Product	RNaseReveal™ Activity Assay Kit, RNaseReveal™ Red Activity Assay Kit
---------	--

Product Description

The RNaseReveal™ Activity Assay Kit is a fluorescent assay for the detection of RNase activity in liquid samples. The buffer and substrate have been formulated for maximum sensitivity, making it a useful tool for detecting RNase contamination in quality control workflows. Biotium also offers [RNase A \(Cat. No. 99871-600uL\)](#) as a separate product that is recommended for use with this kit.

Features

- Exceptional sensitivity, as little as ~0.08 pg RNase
- Simple 40 minute assay
- Detect with handheld fluorometer or fluorescence plate reader
- Useful for detecting RNase contamination in QC workflows

The kit utilizes RNaseReveal™ Substrate which is an RNA probe tagged with a fluorophore and a quencher, so that the intact probe is non-fluorescent. In the presence of RNase, the probe is cleaved and the fluorophore is detached from the quencher, releasing a fluorescent signal. The original RNaseReveal™ Activity Assay has a green fluorescence readout and is more sensitive than RNaseReveal™ Red, with a lower limit of detection of ~ 0.08 pg, using Biotium's [RNase A \(Cat. No. 99871-600uL\)](#) at ≥50 U/mg (Kunitz units). Full substrate cleavage is achieved with ~ 800 pg of the same RNase A. RNaseReveal™ Red features a unique RNase substrate with red fluorescence emission. It was designed for testing samples with intrinsic green fluorescence that interferes with a green fluorescent assay. Both substrates can be detected by any fluorescence plate reader or fluorometer capable of detecting green fluorescence. Biotium also offers [RNase A \(Cat. No. 99871-600uL\)](#) for use with this kit.

Reliable and Sensitive Detection of RNase Contamination

RNaseReveal™ Activity Assay Kit

Kit Name	Abs/Em (nm)	Size	Catalog No.	Kit Components
RNaseReveal™ Activity Assay Kit	496/516	100 assays	31086	RNaseReveal™ Substrate (Cat. No. 31086A)
10X RNaseReveal™ Assay Buffer (Cat. No. 31086B)				
RNaseReveal™ Red Activity Assay Kit	584/604	100 assays	31087	RNaseReveal™ Red Substrate (Cat. No. 31086A)
10X RNaseReveal™ Assay Buffer (Cat. No. 31086B)				

[Learn more about Biotium's unique products for RNA workflows, including RNA extraction kits, quantitation kits, and RNA stains for gel electrophoresis.](#)

This datasheet was generated on January 5, 2026 at 11:34:16 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/rnasereveal-activity-assay-kit/>